

# 3'-O-Me-GTP

## Ribose-modified GTP analogue

3'-O-Me-GTP is an analogue of GTP in which the ribose 3'-hydroxy group has been modified by an ether bond to a methyl group. 3'-O-Me-GTP is an RNA chain-terminating nucleotide that can be useful in DNA transcription studies.

*BLG-M120-25 (5 x 5  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

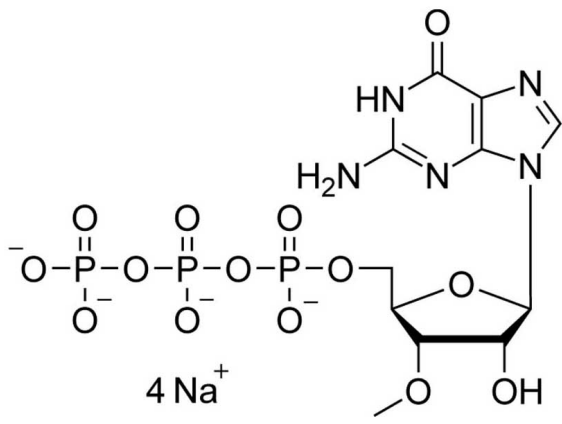
### Ordering Information

[Order Online »](#)

BLG-M120-05	5 $\mu$ mol
BLG-M120-25	5x5 $\mu$ mol

### Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 3'-O-Methylguanosine-5'-O-triphosphate . sodium salt

**Appearance** 10 mM aqueous solution

**CAS** 98923-57-0 (4 Na+) / 61556-45-4 (H+)

**Couple Type** Modified nucleotides

**Excitation Maximum** 252 nm

**Formula**  $C_{11}H_{18}N_5O_{14}P_3$  (free acid)

**MW** 537.2 (free acid)

**Purity** > 95% HPLC

**Technical Info / Product Notes** For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)